

Annexin 5 Protein

Full-length recombinant protein expressed in *E. coli* cells

Catalog # **A61-30H**

Lot # *D135-4*

Product Description

Recombinant full-length human Annexin 5 Protein was expressed by *E. coli* cells using an N-terminal His tag. The gene accession number is [NM_001154](#).

Gene Aliases

ANXA5, PP4, ANX5, ENX2

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 2mM DTT, 25% glycerol.

Storage and Stability

Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

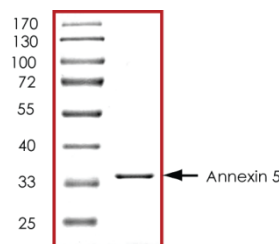
Scientific Background

Annexin 5 is a member of the annexin family of calcium-dependent phospholipid binding proteins. Annexin 5 is a phospholipase A2 and protein kinase C inhibitory protein with calcium channel activity and a potential role in cellular signal transduction, inflammation, growth and differentiation. In vitro studies have shown that Annexin 5 is a cellular protein which plays a role in the inhibition of blood coagulation by competing for phosphatidylserine binding sites with prothrombin (1). Through affinity assays Annexin 5 has been localized on cells that expressed phosphatidylserine on the cell surface, a feature found in apoptosis as well as other forms of cell death (2).

References

1. Koopman, G. et al: Annexin V for flow cytometric detection of phosphatidylserine expression on B cells undergoing apoptosis. *Blood*, 1994; 84 (5): 1415–20.
2. Vermes, I. et al: A novel assay for apoptosis—flow cytometric detection of phosphatidylserine expression on early apoptotic cells using fluorescein labelled Annexin V. *J Immunol Methods*, 1995; 184 (1): 39

Purity



The purity was determined to be **>95%** by densitometry.
Approx. MW **37kDa**.

Annexin 5 Protein

Full-length recombinant protein expressed in *E. coli* cells

Catalog Number	A61-30H
Specific Lot Number	D135-4
Purity	>95%
Concentration	0.2 $\mu\text{g}/\mu\text{l}$
Stability	1 yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.